

File :C:\ISO\Data\23JUN2016\17393-0465774.D
Operator : LauraN
Acquired : 23 Jun 2016 19:50 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-methyl-N,N-Dimethylcathinone (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 7

